

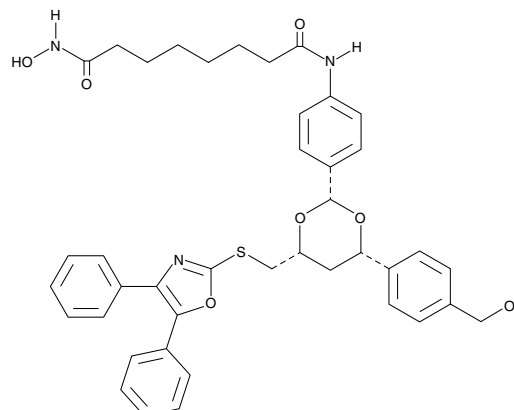
Product Information



Tubacin

Item No. 13691

CAS Registry No.: 537049-40-4
Formal Name: N¹-[4-[(2R,4R,6S)-4-[[[(4,5-diphenyl-2-oxazolyl)thio]methyl]-6-[4-(hydroxymethyl)phenyl]-1,3-dioxan-2-yl]phenyl]-N⁸-hydroxy-octanediamide
MF: C₄₁H₄₃N₃O₇S
FW: 721.9
Purity: ≥90%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 222, 240, 299 nm



Laboratory Procedures

For long term storage, we suggest that tubacin be stored as supplied at -20°C. It should be stable for at least two years.

Tubacin is supplied as a crystalline solid. A stock solution may be made by dissolving the tubacin in the solvent of choice. Tubacin is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of tubacin in these solvents is approximately 16 and 20 mg/ml, respectively.

Tubacin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, tubacin should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Tubacin has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Tubacin is a tubulin acetylation inducer that selectively inhibits histone deacetylase (HDAC) 6 (IC₅₀ = 4 nM).¹ In comparison, it demonstrates at least 300-fold selectivity against all other HDAC isoforms (IC₅₀s range from 1.3-17.3 μM).¹ Tubacin induces α-tubulin hyperacetylation at 2.5 μM in primary cortical neuron cultures.¹ Tubacin inhibition of HDAC6 has been used as a biochemical tool to control microtubule-dependent intracellular trafficking, to manipulate the aggregate formation of misfolded proteins in certain diseases, and to study the dynamics of cellular adhesion.²⁻⁴

References

1. Butler, K.V., Kalin, J., Brochier, C., *et al.* Rational design and simple chemistry yield a superior, neuroprotective HDAC6 inhibitor, Tubastatin A. *J. Am. Chem. Soc.* **132(31)**, 10842-10846 (2010).
2. Chen, S., Owens, G.C., Makarenkova, H., *et al.* HDAC6 regulates mitochondrial transport in hippocampal neurons. *PLoS One* **5(5)**, (2010).
3. Cole, P.A. Chemical probes for histone-modifying enzymes. *Nat. Chem. Biol.* **4(10)**, 590-597 (2008).
4. Xu, K., Dai, X.-L., Huang, H.-C., *et al.* Targeting HDACs: A promising therapy for Alzheimer's disease. *Oxid. Med. Cell Longev.* **2011**, (2011).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/13691

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 07/18/2013

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com